

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Cytokine Panel 1 (mouse) Kit
Lot Number: K0082814
Expiration Date: 31 DEC 2027
Catalog Number: K15245-Series, K152XB-Series, K152WX-Series, K152RD-Series, K152WN-Series, K152NQ-Series, K152QC-Series, K152XC-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (mouse) 96-Well 10-Spot SECTOR Plate	Z0048188	2-8°C	30 APR 2028
Cytokine Panel 1 (mouse) Kit Calibrator	A0080384	2-8°C	28 FEB 2029
Mouse IL-15 Antibody	D0082509	2-8°C	31 MAR 2029
Mouse IL-17A/F Antibody	D0082550	2-8°C	30 APR 2029
Mouse IL-27p28/IL-30 Antibody	D0082549	2-8°C	31 MAY 2029
Mouse IL-33 Antibody	D0082548	2-8°C	30 APR 2029
Mouse IL-9 Antibody	D0082546	2-8°C	30 APR 2029
Mouse IP-10 Antibody	D0082552	2-8°C	30 APR 2029
Mouse MIP-1 α Antibody	D0082551	2-8°C	30 APR 2029
Mouse MIP-2 Antibody	D0082553	2-8°C	30 APR 2029
Mouse/Rat MCP-1 Antibody	D0082547	2-8°C	30 APR 2029
Diluent 41	M0410059	$\leq -10^{\circ}\text{C}$	31 DEC 2027
Diluent 45	M0450056	$\leq -10^{\circ}\text{C}$	29 FEB 2028

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (mouse) Control 1		Cytokine Panel 1 (mouse) Control 2		Cytokine Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-9	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	MCP-1	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	IL-33	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	IL-27p28/IL-30	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
5	IL-15	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	IL-17A/F	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
7	MIP-1 α	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
8	IP-10	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
9	MIP-2	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%



CERTIFICATE OF ANALYSIS

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-9	21.9	75% - 125%	< 20%	2600	75% - 125%	< 20%
2	MCP-1	4.42	75% - 125%	< 20%	325	75% - 125%	< 20%
3	IL-33	1.85	75% - 125%	< 20%	1950	75% - 125%	< 20%
4	IL-27p28/IL-30	5.91	75% - 125%	< 20%	6500	75% - 125%	< 20%
5	IL-15	43.2	75% - 125%	< 20%	26000	75% - 125%	< 20%*
6	IL-17A/F	1.39	75% - 125%	< 20%	1620	75% - 125%	< 20%
7	MIP-1 α	0.378	75% - 125%	< 20%	390	75% - 125%	< 20%
8	IP-10	2.15	75% - 125%	< 20%	650	75% - 125%	< 20%
9	MIP-2	0.577	75% - 125%	< 20%	423	75% - 125%	< 20%

*See Additional Comments

Non-Specific Binding Specifications:

Spot	Name	Spec
1	IL-9	< 1.0%
2	MCP-1	< 1.0%
3	IL-33	< 1.0%
4	IL-27p28/IL-30	< 1.0%
5	IL-15	< 1.0%
6	IL-17A/F	< 1.0%
7	MIP-1 α	< 1.0%
8	IP-10	< 1.0%
9	MIP-2	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-9	3950	pg/mL
2	MCP-1	528	pg/mL
3	IL-33	3010	pg/mL
4	IL-27p28/IL-30	7150	pg/mL
5	IL-15	38000	pg/mL
6	IL-17A/F	2680	pg/mL
7	MIP-1 α	590	pg/mL
8	IP-10	945	pg/mL
9	MIP-2	612	pg/mL

Dilution to top of Curve: 1X

Additional Comments:

Spot	Name	Specification	Comment
5	IL-15	ULOQ CV	Specification for ULOQ is NMT 20%, actual value 24.8% at target of 26000 pg/mL, was passing at 20500 pg/mL.

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 MAR 2026